

Recombinant Human 4-1BB-L

4-1BB-L, also called CD137L or TNFSF9 is a transmembrane glycoprotein belonging to the TNF superfamily (TNFSF) and is involved in reverse signaling that induces apoptosis in T cells and cytokine secretion by monocytes. 4-1BB-L, the ligand for the co-stimulatory immune checkpoint receptor CD137/4-1BB, has been shown to reactivate anergic T lymphocytes in addition to promoting T lymphocyte proliferation. 4-1BB-L through the 4-1BB receptor induces CD4+ and CD8+ T cells to proliferate and secrete cytokines.

Materials Provided

Catalog Number:	92-1321
Quantity Supplied:	1 vial(s), 25 µg per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	32,000 g/mol
Source:	Recombinant
Purity:	>97%
Endotoxin Level:	<0.10 EU
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 µg/mL
Reconstitution Volume:	0.250 mL
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com